

Identification of peptide L, a novel neuropeptide that regulates the expression of L-type voltage-gated calcium channels in photoreceptors

174.03/LL2



Liheng Shi¹, Michael Ko¹, Louise Abbott¹, Gladys Y.-P.Ko¹

¹ Department of Veterinary Integrative Biosciences, College of Veterinary Medicine and Biomedical Sciences, Texas A&M University, College Station, TX 77843-4458, USA. (Gladys Ko: gko@cvm.tamu.edu)



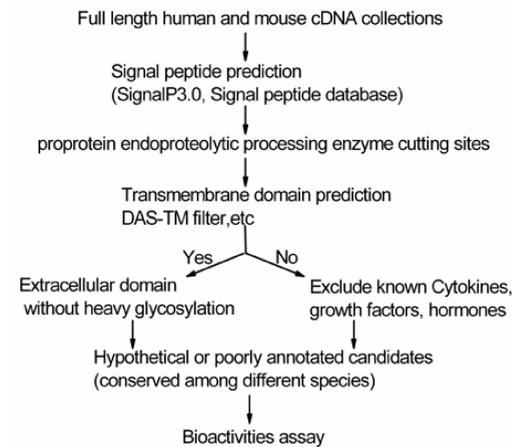
Introduction:

- Neuropeptides act like peptidic hormones or neurotransmitters and play diverse roles in regulating neural functions (1).
- Computational bioinformatics has become a powerful tool to discover novel bioactive peptides (2).
- The L-type voltage-gated calcium channels (L-VGCCs) mediate a voltage-dependent, depolarization-induced Ca²⁺ influx and regulate diverse biological processes such as contraction, secretion, differentiation, synaptic plasticity, and gene expression (3).
- The L-VGCCs are essential in gating pro-longed neurotransmitter release in retinal photoreceptors (4).

Methods:

Cloning of mouse peptide L full length cDNA
Cell culture and transfection
High-performance liquid chromatography (HPLC) and mass spectrometry (MS)
In situ hybridization
Immunoblotting, quantitative (Q)-PCR, and cAMP enzyme immunoassay
GST fusion protein purification and peptide synthesis
Patch Clamp Electrophysiology

Figure 1. A flow chart illustration of our *in silico* computational screening strategy and procedure.



Human and mouse full length protein databases were subjected to a separation with five steps by computer algorithms including N-terminal signal peptide sequence (secretion marker), propeptide convertase targeting motif, transmembrane domain, potential glycosylation modification, and sequence homology across species. Potential candidates were ranked according to the score assigned from each step.L (5805.6982 m/z). The 49 amino acid sequence of peptide L is listed at the top of this panel.

Figure 2. Sequence alignment between human and mouse propeptide L.

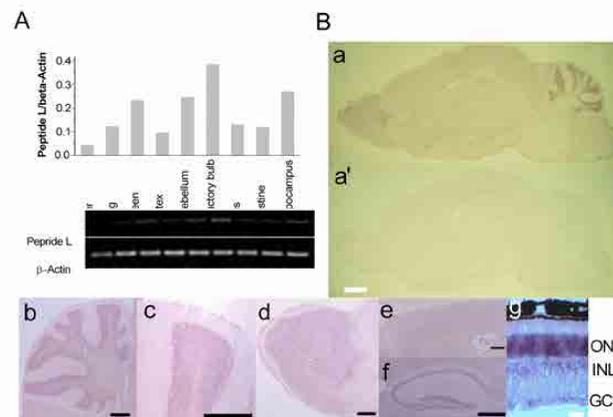
Protein sequence alignment

Human	--MRLLLAALAAALARAPAEVCAALNVTVSPGPPVDYLEGENATLLCHVSQKRRKDSLL	58
Mouse	MRLRLLLAALAAVLLG--PAPEVCGALNVTVSPGPPVDYLEGENATLLCHVSQKRRKDSLL	58
Human	AVIWFYFAPSFDSCQKLLWVNTLRLVQVYGHFSSRSARRRLLRLLERQALYRLSVLTLQ	118
Mouse	AVIWFYFAPD-DSQKLLWVNTLRLVQVYGHFSSRSARRRLLRLLERQALYRLSVLTLQ	117
Human	PSDQGHYVCRVQEISRHRNKWTAWNGSSATEMRVLSLKASESSPEKTKETWAFEDLLY	178
Mouse	PTDQGHYVCRVQEISRHRNKWTAWNGSSATEMRVLSLKASESSPEKTKETWAFEDLLY	177
Human	VYAVLVCCVGLLSILLFMLVYVWQSVFNRKRSRVRHYLVKCPQNSGGETVTSVTSLAPLQ	238
Mouse	VYAVLVCCVGLLSVLLFPLVIAWQSVFNRKRSRVRHYLVKCPQNSGGETVTSVTSLAPLQ	237
Human	PKGKRQKPKDIPPAVPAKAPIATPFHPKPLLPQRKVTLPRIASENLYAELELKP	298
Mouse	PQKGRQKPKDIPPAVPAKAPIATPFHPKPLLPQRKVTLPRIASENLYAELELKP	297
Human	RAAKGAPTSTVYAQLFEENKL	320
Mouse	RAAKGAPTSTVYAQLFEENQL	319

■ Signal peptide ■ Propeptide convertase cutting site
■ Transmembrane domain ■ Peptide L ■ Peptide II

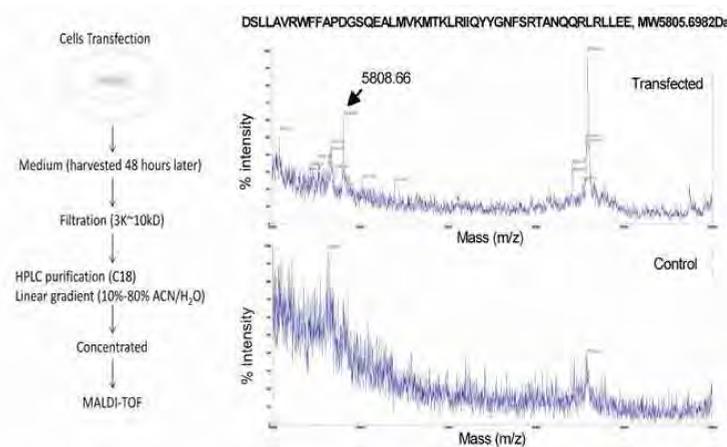
Human and mouse peptide L proprotein are over 80% homologous. The sequence contains a signal peptide sequence (green), proprotein convertase cutting motifs (red), the main peptide L coding region (blue), and a transmembrane domain (purple). Peptide II (control for the electrophysiological studies) is highlighted in orange.

Figure 3. Expression of peptide L in mouse tissues.



(A) Gene expression of peptide L in various mouse tissues was determined by RT-PCR. The mRNA of peptide L was detected in the liver, lung, spleen, intestine, eyes, cerebral cortex (cortex), cerebellum, olfactory bulb, and hippocampus, with relatively higher expression levels in the spleen, cerebellum, olfactory bulb, and hippocampus. (B) The mRNA expression of peptide L in the mouse brain and retina was detected by *in situ* hybridization. Adult mouse brain sagittal sections and retina sections were cut at 10 μ m thickness and processed for *in situ* hybridization. The sense probe served as the negative control (a'). Peptide L is strongly expressed in the olfactory bulb (a, d), cerebellum (a, d, c), cerebral cortex (a, e), and hippocampus (a, c). In the mouse retina (g), the expression of peptide L was detected in the outer nuclear layer (ONL, strongest signal), inner nuclear layer (INL), and ganglion cell layer (GCL). The scale bars represent 1.0 mm (a and a'), 0.1 mm (b, c, d, e, f), and 0.02 mm (g).

Figure 4. Verification of peptide L secretion by MALDI-TOF mass spectrometry.

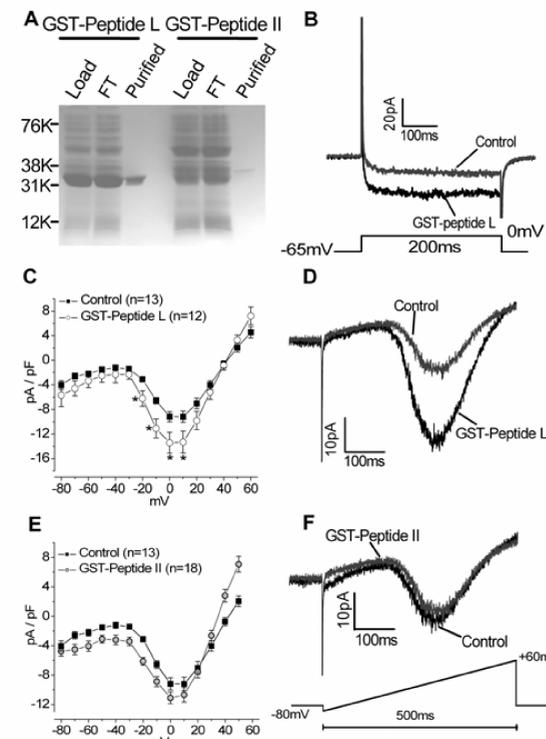


Mouse 661w cells were transfected with the propeptide L gene (mouse E130203B14Rik) expression vector or empty vector (control). Culture media were collected 36-40 hrs after transfection, filtered (3~10kD), and subjected to HPLC. The final elution (linear gradient 10%-80% ACN/H₂O and 0.1% TFA in all solutions) was concentrated by speedvac and analyzed by MALDI-TOF MS. The arrowhead indicates the peak (5808.66 m/z) that corresponds to the predicted molecular weight of peptide L (5805.6982 m/z). The 49 amino acid sequence of peptide L is listed at the top of this panel.

Acknowledgements:

This work was supported by NIHRO1 EY017452 to GK and a postdoctoral research award from the College of Veterinary Medicine and Biomedical Sciences at Texas A&M University to LS. We thank Drs. Yohannes Rezenom and William Russell, Laboratory for Biological Mass Spectrometry, Texas A&M University, for their technical support.

Figure 5. GST-peptide L increased L-VGCC currents in cone photoreceptors.



GST-peptide L and GST-peptide II (control) were expressed and purified from *E. coli* (A). Cultured E18 cone photoreceptors were treated with PBS (control), 800 ng/ml GST-peptide II (control peptide), or 800 ng/ml GST-peptide L for 4 hr prior to recordings. Whole cell patch clamp recordings were performed under both ramp and step commands (B, D). Representative current traces under a ramp command (D, F) or a step command (E) are shown. Average current-voltage (I-V) relationships were obtained from the step command as in current density (pA/pF), and the maximum inward current densities were elicited at 0-10 mV of the step command (C, E). The average maximal current density obtained from GST-peptide L treated cells (n=12; open circle, C) was significantly larger than the control (n=13; filled square). No significant difference occurred between the control and GST-peptide II treated cells (n=18; gray circle; E). Comparisons between the control and recombinant protein groups were made using Student's *t*-test; **p* < 0.05.

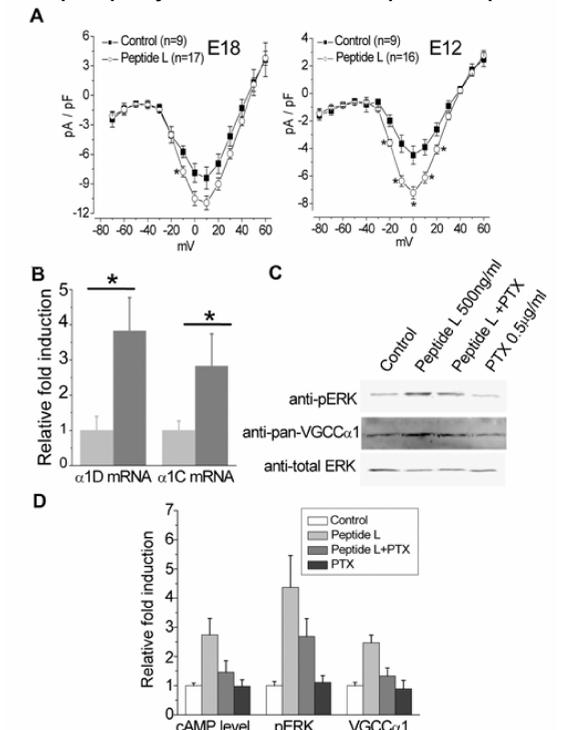
Conclusion:

- Computational bioinformatics coupled with electrophysiological activities assay can be an effective method to identify bioactive neuropeptides.
- Peptide L up-regulated the expression of L-VGCCs in photoreceptors.
- Peptide L might be important in modulating neurotransmitter release in photoreceptors.

References:

1. Rotzinger, S., Lovejoy, D. A., and Tan, L. A. (2010) *Peptides* 31, 736-756
2. Mirabeau, O., Perlas, E., Severini, C., Audero, E., Gascuel, O., Possenti, R., Birney, E., Rosenthal, N., and Gross, C. (2007) *Genome Res* 17, 320-327
3. Catterall, W. A., and Few, A. P. (2008) *Neuron* 59, 882-901
4. Barnes, S., and Kelly, M. E. (2002) *Adv Exp Med Biol* 514, 465-476
5. Ko, M. L., Liu, Y., Dryer, S. E., and Ko, G. Y. (2007) *J Neurochem* 103, 784-792

Figure 6. Synthesized peptide L enhanced the protein expression L-VGCC α 1 subunits, cAMP production, and ERK phosphorylation in chicken cone photoreceptors.



Cultured chick photoreceptors (E10+2 and E16+2) were treated with 500 ng/ml synthesized peptide L or control buffer for 4 hrs prior to patch clamp recordings (A). Averaged current-voltage (I-V) relationships were obtained from the step command as in current density (pA/pF), and the maximum inward current densities were elicited at 0-10 mV of the step command. Peptide L increased current density at both E12 and E18 (compared to their respective control groups). E12: control, -4.49 \pm 0.66 pA/pF (n=9, dark square); peptide L treated cells, -7.22 \pm 0.45 pA/pF (n=16, open square). E18: control, -8.40 \pm 1.11 pA/pF (n=9, dark square); peptide L, -10.93 \pm 0.70 pA/pF (n=17, open square). The mRNA levels of both L-VGCC α 1C and α 1D were measured by Q-PCR (B; n=6 for each group). Comparisons between the control and peptide L groups were made using Student's *t*-test; **p* < 0.05. L-VGCC α 1, ERK, and pERK were detected by western blot. The band densities were measured and relative values were calculated as a ratio to total ERK (C). Relative fold difference (n=3) was calculated by comparison to control samples which were set to one (D). Peptide L significantly increased cAMP levels, phosphorylation of ERK, and L-VGCC α 1 expression (D, gray bar). However, PTX, a specific GPCR inhibitor, blocked peptide L effects (D, dark gray bar). Values are means \pm sem.